



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: CHAPS 5% Solution (Filtered)

Catalog No: DG098

Release Date: 1/29/20

Size: 100ml

Retest Date: 2/29/25

Lot No: 202901

Storage: 4°C

Specifications

CAS Number: N/A

Appearance: Clear, colorless solution

Test	Specification	Result
Conductivity	2.2ppm / 4.38μS	PASS
pH	4.6 – 5.8	5.42

This document has been produced electronically and is valid without signature

Date of Print: February 18, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com